



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## Renilla Luciferase Assay Kit 2.0

Highly sensitive & linear flash-type luminescence assay for measuring *Renilla* luciferase activity in transfected cell reporter gene assays.



### Product attributes

<b>Assay type/options</b>	Endpoint assay, Luminescence (flash-type)
<b>Detection method/readout</b>	Luminometer (single-tube or microplate reader with reagent injectors)
<b>Storage Conditions</b>	Store at -60 °C or below

## Product Description

This kit is designed to measure *Renilla* luciferase (RLuc) activity in transfected cell reporter gene assays.

*Renilla* luciferase has been used as a reporter gene for studying gene regulation and function *in vitro* and *in vivo*. It commonly is used in multiplex transcriptional reporter assays or as a normalizing transfection control for firefly luciferase assays. *Renilla* luciferase, a monomeric 36,000 Dalton protein, catalyzes coelenterazine oxidation by oxygen to produce light. The enzyme does not require post-translational modification for its activity and may function as a genetic reporter immediately following translation.

Coelenterazine also emits light from enzyme-independent oxidation, a process known as autoluminescence. The autoluminescence is enhanced by superoxide anion and peroxynitrite in cells and tissues. This assay kit utilizes a special buffer formulation designed to yield reliable, linear measurements of *Renilla* luciferase activity with minimal autoluminescence background and superior sensitivity.

**Note:** This kit is a flash-type luminescence assay, which requires luminescence to be measured immediately after adding reagents to your sample. It requires a single-tube luminometer or a luminescence microplate reader with reagent injectors.

### Easy-to-Use for Consistent Results

This kit is a new and improved replacement for the *Renilla* Luciferase Assay Kit (catalog no. 30004). In addition to providing a simplified protocol, the *Renilla* Luciferase Assay Kit 2.0 features Biotium's Aquaphile™ Coelenterazine, a water soluble substrate that can be stored at -20 °C with minimal evaporation, unlike methanol solutions of coelenterazine.

The stated number of assays is based on 96-well assay format. The kits include enough Passive Lysis Buffer to perform the stated number of assays in plate sizes ranging from 96-well to 24-well. If needed, additional 5X Passive Lysis Buffer (catalog no. 99912) can be purchased separately (see below).

Biotium also offers the [Firefly & \*Renilla\* Luciferase Single Tube Assay Kit](#) (catalog no. 30081), a combined luciferase assay kit for sequential measurement of Firefly and *Renilla* luciferase activity sequentially in the same sample with high sensitivity and linearity.

## More Luciferase Kits & Related Products

Biotium Cat. No.	Product	Description	Promega Assay	Promega Cat. No.
<a href="#">30138</a>	<a href="#">Steady-ATP™ HTS Viability Assay Kit</a>	Highly sensitive ATP-based cell luminescent viability assay with signal half-life of 5 hours for high-throughput screening in multi-well plates.	CellTiter-Glo® Luminescent Cell Viability Assay	G7570, G7571, G7572, G7573
<a href="#">30085</a>	<a href="#">Firefly Luciferase Assay Kit 2.0</a>	Flash-type firefly luciferase assay kit.	Firefly Luciferase Assay System	E1483, E1500, E1501, E4030, E4530, E4550
<a href="#">30075</a>	<a href="#">Firefly Luciferase Assay Kit (Lyophilized)</a>	Firefly luciferase assay kit with lyophilized assay buffer for economical room temperature shipping and convenient storage at -20°C.		
<a href="#">30028</a>	<a href="#">Steady-Luc&amp;#153; Firefly HTS Assay Kit</a>	Glow-type firefly luciferase assay with signal half-life of about 3 hours for high-throughput screening in multi-well plates.	Steady-Glo® Assay System	E2510, E2520, E2550
<a href="#">30028L</a>	<a href="#">Steady-Luc Firefly HTS Assay Kit (Lyophilized)</a>	Steady-Luc kit with lyophilized assay buffer for economical room temperature shipping and convenient storage at -20°C.		
<a href="#">30082</a>	<a href="#">Renilla Luciferase Assay Kit 2.0</a>	Flash-type Renilla luciferase assay kit.	<i>Renilla</i> Luciferase Assay System	E2810, E2820
<a href="#">30081</a>	<a href="#">Firefly and Renilla Luciferase Single Tube Assay Kit</a>	Assay kit for measuring firefly and Renilla activities sequentially in the same sample in a single tube.	Dual-Luciferase® Assay System (Firefly & <i>Renilla</i> )	E1910, E1960, E1980

This datasheet was generated on June 7, 2026 at 06:05:47 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/renilla-luciferase-assay-kit-2-0/>